



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-04623

Product Description: Cytokine Panel 1 (human) Control Pack

Catalog Number: C4050-1

Lot Number: CP000018

Expiration Date: 31 MAY 2019

Concentration of Materials Tested:

Units = pg/mL	Assay									
	hu GM-CSF	hu IL-1 α	hu IL-5	hu IL-7	hu IL-12/23 p40	hu IL-15	hu IL-16	hu IL-17A	hu TNF- β	hu VEGF
Reconstituted Control 1 (A00C0251)	285	135	303	232	1122	191	1208	1760	168	338
Reconstituted Control 2 (A00C0252)	67.1	30.2	76.5	58.5	257	48.3	261	431	40.4	72.3
Reconstituted Control 3 (A00C0263)	21.4	13.7	64.1	17.8	84.2	16.2	148	210	11.5	50.2

Reconstitution Volume: 250 μ L

Control concentrations were assigned across a minimum of three kit lots.

Cytokine Panel 1 (human) Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 1 (human) Control 1	A00C0251	2-8°C	31 MAY 2019
Cytokine Panel 1 (human) Control 2	A00C0252	2-8°C	31 MAY 2019
Cytokine Panel 1 (human) Control 3	A00C0263	2-8°C	31 JUL 2019

Statement: The above product was manufactured and tested using MSD procedures and it conforms to MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Caitlin Hansen	Quality Control		25 JUL 2014